



Auto-Pure 20B

Automated cell isolation system

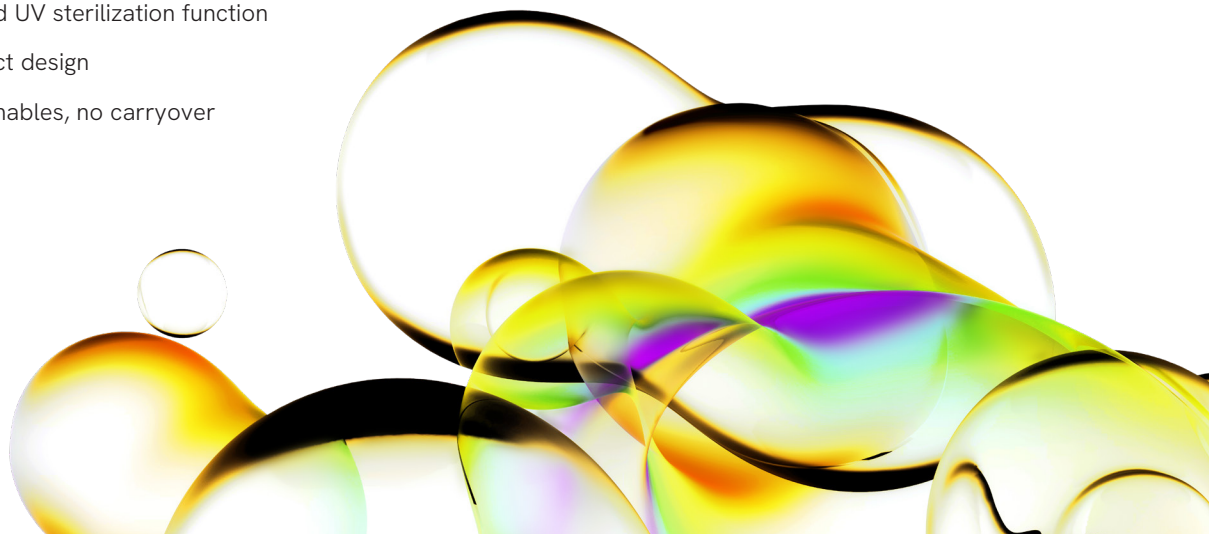
The Auto-Pure 20B serves as a versatile automated cell isolation platform for the T-SPOT.TB test. It's an ideal solution for laboratories with low throughput requirements, while also providing valuable backup support in high-volume testing environments.



A validated protocol developed by Revvity enables users to quickly and efficiently automate the important peripheral blood mononuclear cell (PBMC) isolation stage of the **T-SPOT™.TB** test. The Auto-Pure 20B uses magnetic rods to transport PBMCs through the various purification phases; mixing, binding, washing and elution.

Built in flexibility

The Auto-Pure 20B cell isolation system utilises a strip based workflow, bringing enhanced operational efficiency to PBMC isolation. This allows users to deal with runs of variable sizes, optimising consumable usage.

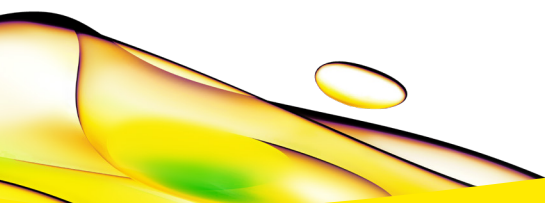
- Easy to use with minimal maintenance
- Simple touch screen interface
- Efficient hands-free processing: handles up to 20 samples in under 45 minutes
- Drawer design for easy loading
- Pre-loaded validated protocol
- Internal lighting and UV sterilization function
- Bench-top, compact design
- Disposable consumables, no carryover or contamination



Model	Auto-Pure 20B	
Features	AUTOPURE 20B	
Processing volume	3.5 mL*	
Processing time	Less than 45 minutes	
Number of samples	1-20	
Protocol storage	Pre-loaded cell isolation protocol	
External interfaces	Standard USB, Ethernet port, Wi-Fi, Built-in SD card	
Operating conditions	Ambient temperature: 10 - 35 °C Relative humidity: ≤ 70 % Input voltage: AC 100 - 240 V, 50 Hz/60 Hz	
Service information	2 levels of service contract available (Gold/Silver)	
Power supply	450 W	
Dimensions (WxDxH)	400 × 520 × 450 mm	
Weight	30 kg	
Regulatory status	CE IVD and FDA approved	
Consumables	Order code	
AP20B/10B_Tube Strips	TUBESTRIPS10/20B	
AP20B/10B_Magnetic Rod Tips	AUTOPURETIPCOMB	

*Could vary depending on age and immune status

Products may not be licensed in accordance with the laws in all countries.
Please check with your local representative for availability.



revvity